

PRODUCT INFORMATION



JNJ-64619178

Item No. 29231

CAS Registry No.: 2086772-26-9

Formal Name: (1S,2R,3S,5R)-3-[2-(2-amino-3-bromo-7-quinoliny)ethyl]-5-(4-amino-7H-pyrrolo[2,3-d]pyrimidin-7-yl)-1,2-cyclopentenediol

MF: C₂₂H₂₃BrN₆O₂

FW: 483.4

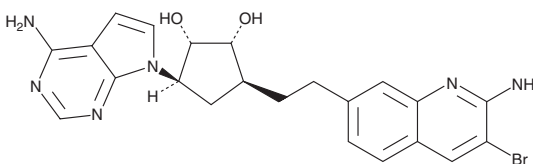
Purity: ≥98%

UV/Vis.: λ_{max}: 214, 233 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

JNJ-64619178 is supplied as a solid. A stock solution may be made by dissolving the JNJ-64619178 in the solvent of choice, which should be purged with an inert gas. JNJ-64619178 is soluble in DMSO.

Description

JNJ-64619178 is an inhibitor of protein arginine methyltransferase 5 (PRMT5; IC₅₀ = 0.14 nM).¹ It inhibits the growth of various cancer cells *in vitro* and reduces tumor growth in non-small cell lung cancer (NSCLC) and small cell lung cancer (SCLC) mouse xenograft models.²

References

1. Tao, H., Yan, X., Zhu, K., *et al.* Discovery of novel PRMT5 inhibitors by virtual screening and biological evaluations. *Chem. Pharm. Bull. (Tokyo)* **67**(4), 382-388 (2019).
2. Wu, T., Millar, H., Gaffney, D., *et al.* Abstract 4859: JNJ-64619178, a selective and pseudo-irreversible PRMT5 inhibitor with potent *in vitro* and *in vivo* activity, demonstrated in several lung cancer models. *Cancer Res.* **78**(13_suppl.), 4859 (2018).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/10/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM